



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** Alamar Blue Cell Viability Reagent

**Catalog #:** 786-921

**Lot #:** 190212

**Re-test date:** 02-2024

### Tests:

Appearance

Dark blue solution with no visible ppt.

Release Method

Absorbance 0.74-0.83 AU

Microbiology testing

No growth at 34°C to 36°C for minimum of 72 hours

**Storage:** Store at 4°C

*for Research Use Only*

This document has been produced electronically and is valid without signature

**Date of Print:** March 24, 2020

### **G-Biosciences / Genotech**

9800 Page Avenue

St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504

Alamar blue/Com#3  
Rev 04.14.17CCH



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)